

## Batch information

### control lysate

HTRF MLKL Total Kit - Ctrl lysate

Product reference :

**64MLKLTTDA**

### Product characteristics :

- Lot Number	01A
- Conditioning buffer	Lysis buffer 4 + blocking reagent
- Appearance	Frozen form
- Storage conditions	-60 °C
- Expiration date	2024.07.12

### Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ab)	150.00 µl	N/A

**Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: [www.cisbio.com](http://www.cisbio.com)**

